

---

# DCAMKL1

Product code: 02-139

Full-length human DCAMKL1 [1-729(end) amino acids of accession number NP\_004725.1] was expressed as N-terminal GST-fusion protein (108 kDa) using baculovirus expression system.

Assay platform : MSA  
Substrate : GS peptide

ATP ( $\mu$ M) Km app / Bin : 230 / 250  
Metal : Mg  
Reference compound : Staurosporine  
IC50 at ATP Bin (nM) : 200  
IC50 at 1 mM ATP (nM) : 410

---